



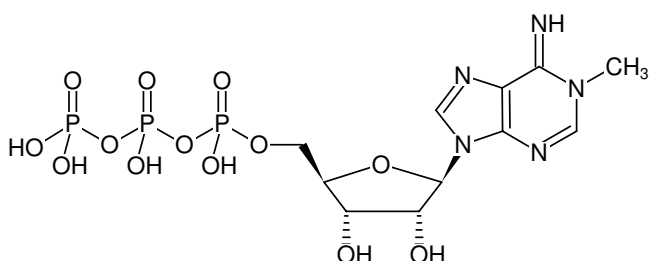
N¹-Methyl-ATP - Solution

N¹-Methyl-adenosine-5'-triphosphate, Sodium salt

Cat. No.	Amount
NU-987S	10 µl (100 mM)
NU-987L	5 x 10 µl (100 mM)

Selected References:

Dominissini *et al.* (2016) The dynamic N1-methyladenosine methylome in eukaryotic messenger RNA. *Nature* **530** (7591):441.



Structural formula of N¹-Methyl-ATP - Solution

For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Short term exposure (up to 1 week cumulative) to ambient temperature possible.

Shelf Life: 12 months after date of delivery

Molecular Formula: C₁₁H₁₈N₅O₁₃P₃ (free acid)

Molecular Weight: 521.21 g/mol (free acid)

Exact Mass: 521.01 g/mol (free acid)

Purity: ≥ 95 % (HPLC)

Form: solution in water

Color: colorless to slightly yellow

Concentration: 100 mM - 110 mM

pH: 7.5 ±0.5

Spectroscopic Properties: λ_{max} 257 nm, ε 14.6 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)